

Rat MIP-1 alpha ELISA Kit (CCL3) ab213916

[1 References](#) [1 Image](#)

Overview

Product name Rat MIP-1 alpha ELISA Kit (CCL3)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	18pg/ml	1.08	= 6%
2	16	54pg/ml	3.61	= 6.7%
3	16	267pg/ml	15.75	= 5.9%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	18pg/ml	1.11	= 6.2%
2	24	56pg/ml	4.48	= 8%
3	24	248pg/ml	14.88	= 6%

Sample type

Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

Assay type

Sandwich (quantitative)

Sensitivity

< 1 pg/ml

Range

7.8 pg/ml - 500 pg/ml

Assay time

3h 30m

Assay duration

Multiple steps standard assay

Species reactivity

Reacts with: Rat

Product overview

The MIP-1alpha Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213916) is designed for the quantitative measurement of MIP-1 alpha in cell culture supernatants, serum and plasma (heparin, EDTA, citrate).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for MIP-1 alpha has been pre-coated onto 96-well plates. Standards and test samples are added to the wells; a biotinylated detection polyclonal antibody from goat specific for MIP-1 alpha is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the MIP-1 alpha amount of sample captured in plate.

Notes Macrophage inflammatory protein-1 alpha (MIP-1 alpha), also called CCL3, LD78. The cDNA for MIP-1 alpha predicts a mature peptide of 69 amino acids with a molecular mass of 7,889 daltons. LD78 is a member of a newly identified superfamily of small inducible proteins involved in inflammatory responses, wound healing and tumorigenesis. MIP-1 alpha is a chemokine that has pro-inflammatory and stem cell inhibitory activities in vitro. It constitutes an important second signal for mast cell degranulation in the conjunctiva in vivo and consequently for acute-phase disease.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-rat MIP-1 alpha coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- rat MIP-1 alpha antibody		1 x 100µl
Lyophilized recombinant rat MIP-1 alpha standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

Function Monokine with inflammatory and chemokinetic properties. Binds to CCR1, CCR4 and CCR5. One of the major HIV-suppressive factors produced by CD8+ T-cells. Recombinant MIP-1-alpha induces a dose-dependent inhibition of different strains of HIV-1, HIV-2, and simian immunodeficiency virus (SIV).

Sequence similarities Belongs to the intercrine beta (chemokine CC) family.

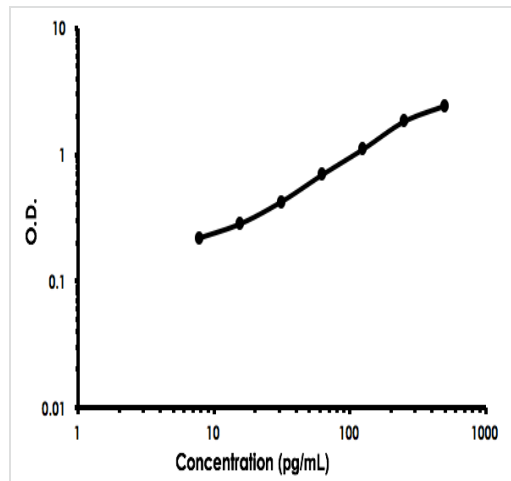
Post-translational modifications

N-terminal processed form LD78-alpha(4-69) is produced by proteolytic cleavage after secretion from HTLV1-transformed T-cells.

Cellular localization

Secreted.

Images



MIP-1 alpha ELISA Kit (ab213916) Standard Curve

MIP-1 alpha ELISA Kit (ab213916) Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors